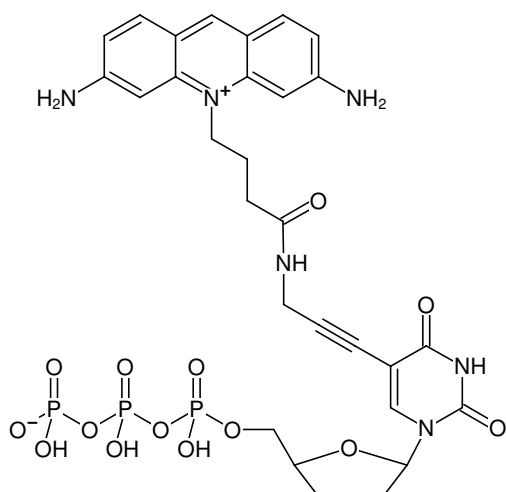




### 5-Propargylamino-ddUTP-ATTO-465

5-Propargylamino-2',3'-dideoxyuridine-5'-triphosphate, labeled with ATTO 465, Triethylammonium salt

Cat. No.	Amount
NU-1619-465	160 µl (0,5 mM)



Structural formula of 5-Propargylamino-ddUTP-ATTO-465

**For general laboratory use.**

**Shipping:** shipped on gel packs

**Storage Conditions:** store at -20 °C

Short term exposure (up to 1 week cumulative) to ambient temperature possible.

**Shelf Life:** 12 months after date of delivery

**Molecular Formula:** C<sub>29</sub>H<sub>33</sub>N<sub>6</sub>O<sub>14</sub>P<sub>3</sub> (free acid)

**Molecular Weight:** 782.53 g/mol (free acid)

**Exact Mass:** 782.13 g/mol (free acid)

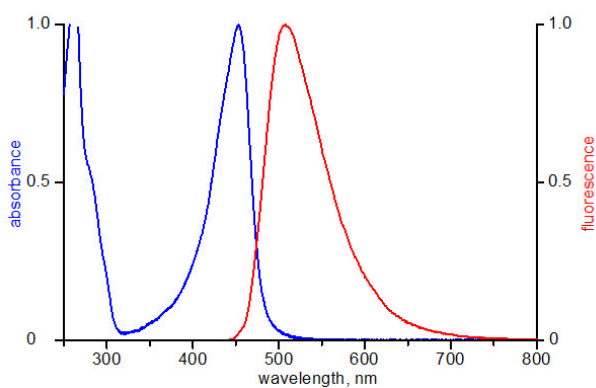
**Purity:** ≥ 95 % (HPLC)

**Form:** solution in water

**Concentration:** 0.50 mM - 0.55 mM

**pH:** 7.5 ± 0.5

**Spectroscopic Properties:** λ<sub>exc</sub> 453 nm, λ<sub>em</sub> 506 nm, ε 75.0 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5)



excitation and emission spectrum of ATTO 465